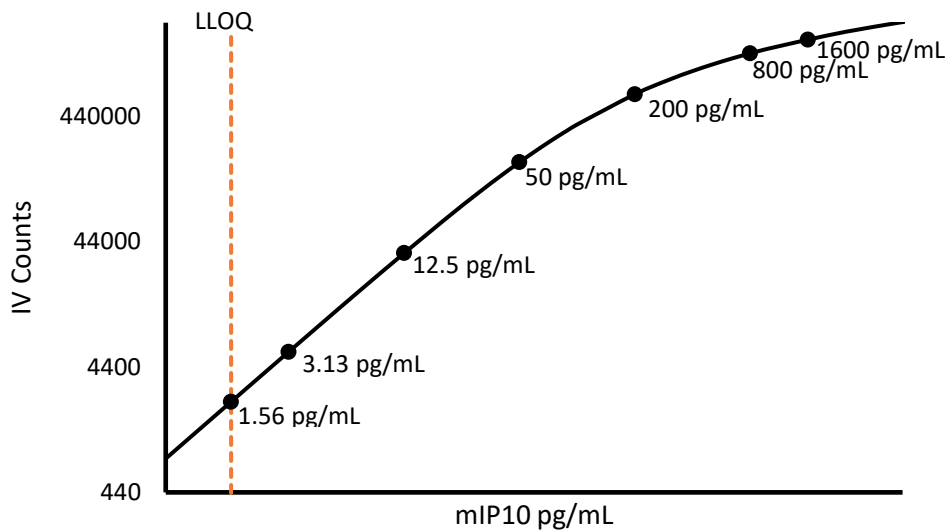


Description – Mouse IP-10

IFN-γ-inducible protein 10 (IP-10, CXCL10) is a 10 kDa chemokine secreted from cells stimulated with type I and II interferons (IFNs) and lipopolysaccharide (LPS). IP-10 is constitutively expressed at low levels in thymic, splenic, and lymph node stroma. Elevated levels of IP-10 protein have been found in the cerebral spinal fluid in patients with viral meningitis and multiple sclerosis. In these diseases, levels of IP-10 correlate with the tissue infiltration of T lymphocytes, suggesting that IP-10 plays an important role in the recruitment of T cells to sites of tissue inflammation.

Calibration Curve: Calibrator concentrations and Lower Limit of Quantification are depicted in the figure below. This standard curve is for demonstration purposes; end users should prepare a standard curve for each assay run.



Minimum Required Dilution (MRD)

Diluted Sample volume (1:2 Dilution)*	50 µL per measurement
--	-----------------------

*See Kit Instructions for details

Endogenous Serum and Plasma Readings: Healthy EDTA plasma and serum samples (n=8) from non-medicated, non-immunized mice were measured.

% Above LOD	100%
% Above LLOQ	100%

Assay Range: The upper end of the dynamic range is equal to the top calibrator concentration multiplied by MRD.

Analytical LLOQ	1.56 pg/mL
Functional LLOQ (x MRD)	3.12 pg/mL
LOD	0.618 pg/mL
Assay Range	0 – 3200 pg/mL

Note: Data described were developed during assay development. Under different assay conditions, assay may perform differently than shown. For complex matrices such as serum or plasma, assay diluent optimization (for example by adding blocking agents) may improve performance of these matrices in this assay.